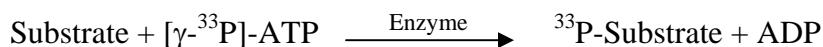


## JAK2 (V617F)

(Tyrosine-protein kinase JAK2, Janus kinase 2, JAK-2)

- **CAT#:** JAK2 (V617F)
- **Enzyme:** Human JAK2 (V617F)
- **Substrate:** Peptide substrate, poly[Glu:Tyr] (4:1), 0.2 mg/ml  
ATP 10  $\mu$ M
- **Reaction:**



- **Reference Compound IC50s:**

Compounds	IC50 (nM)
Staurosporine	<0.1
D-64406	27,000

- **Scientific Information:** O60674 (<http://www.expasy.org/uniprot/O60674>)

### Reference compound IC50 for JAK2 (V617F)

